



6042 Cornerstone Court West, Suite B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

**ATP (500  $\mu$ M)**  
Catalog #: 79686

**Description:** Suitable for Kinase-Glo and ADP-Glo detection and for use with select BPS Bioscience Kinase Assay Kits.

**Supplied as:** Aqueous buffer solution.

**Concentration:** Final concentration should be 10-20  $\mu$ M in assay.

**Warning:** Avoid freeze/thaw cycles.

**Size:** 200  $\mu$ l

**Storage:** -20°C

### Optimized for use with:

Catalog #	Name	Size
79576	<a href="#">ABL (T315I) Assay Kit</a>	96 reactions
79549	<a href="#">ALK1 Kinase Assay Kit</a>	96 reactions
79605	<a href="#">ALK2 Kinase Assay Kit</a>	96 reactions
48688	<a href="#">B-Raf(V600E) Kinase Assay Kit</a>	96 reactions
79568	<a href="#">BTK Assay Kit</a>	96 reactions
79660	<a href="#">c-Mer Kinase Assay Kit</a>	96 reactions
79559	<a href="#">c-Met Kinase Assay Kit</a>	96 reactions
79570	<a href="#">c-Raf Kinase Assay Kit</a>	96 reactions
79597	<a href="#">CDK1 Assay Kit</a>	96 reactions
79599	<a href="#">CDK2 Assay Kit</a>	96 reactions
79674	<a href="#">CDK4 Assay Kit</a>	96 reactions
79600	<a href="#">CDK5 Assay Kit</a>	96 reactions
79603	<a href="#">CDK7 Assay Kit</a>	96 reactions
79628	<a href="#">CDK9 Assay Kit</a>	96 reactions

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court West, Suite B  
 San Diego, CA 92121  
 Tel: 1.858.202.1401  
 Fax: 1.858.481.8694  
 Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

40326	<a href="#">EGFR (T790M/C797S/ L858R) Kinase Assay Kit</a>	96 reactions	
40321	<a href="#">EGFR Kinase Assay Kit</a>	96 reactions	
40324	<a href="#">EGFR(L858R) Kinase Assay Kit</a>	96 reactions	
40323	<a href="#">EGFR(T790M) Kinase Assay Kit</a>	96 reactions	
40322	<a href="#">EGFR(T790M/L858R) Kinase Assay Kit</a>	96 reactions	
40722	<a href="#">FAK Kinase Assay Kit</a>	96 reactions	
40721	<a href="#">HER2 Kinase Assay Kit</a>	96 reactions	
79518	<a href="#">JAK1 (Janus Kinase 1) Assay Kit</a>	96 reactions	
79520	<a href="#">JAK2 (Janus Kinase 2) Assay Kit</a>	96 reactions	
79521	<a href="#">JAK3 (Janus Kinase 3) Assay Kit</a>	96 reactions	
79566	<a href="#">RET Assay Kit</a>	96 reactions	
79560	<a href="#">RIPK1 Kinase Assay Kit</a>	96 reactions	
79689	<a href="#">RIPK1 Kinase Assay Kit (384-well)</a>	384 reactions	
79680	<a href="#">SRC Assay Kit</a>	96 reactions	
79671	<a href="#">SYK Assay Kit</a>	96 reactions	
79548	<a href="#">TrkA Assay Kit</a>	96 reactions	
79527	<a href="#">TYK2 (Tyrosine Kinase 2) Assay Kit</a>	96 reactions	
79502	<a href="#">TYR Phosphorylation Assay Kit</a>	96 reactions	
79593	<a href="#">TYRO3 Kinase Assay Kit</a>	96 reactions	
40325	<a href="#">VEGFR2 (KDR) Kinase Assay Kit</a>	96 reactions	
79681	<a href="#">YES1 Assay Kit</a>	96 reactions	

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.  
 To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)